



CERTIFICATE OF ANALYSIS

Product Description: S-PLEX Human Tau (pT217) Kit
Lot Number: K00S0276
Expiration Date: 30 APR 2027
Catalog Number: K151APFS-Series; K156APFS-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Biotin Human Tau (pT217) Antibody	B00T0257	2-8°C	31 AUG 2028
TURBO-BOOST Human Tau Antibody	D00T0257	2-8°C	31 OCT 2028
Blocker S1	R0140016	≤ -10°C	31 DEC 2027
Diluent 2	M0020133	≤ -10°C	30 JUN 2027
S-PLEX Enhance E1	Y0550024	≤ -10°C	31 OCT 2027
S-PLEX Enhance E2	Y0560024	≤ -10°C	31 DEC 2027
Human Tau (pT217) Calibrator	A00S0121	≤ -70°C	30 APR 2027
S-PLEX Coating Reagent C1	521330	≤ -70°C	30 JUN 2028
S-PLEX Detect D1 with TURBO-TAG™ Reagent	Y0580034	≤ -70°C	31 DEC 2027
S-PLEX Detect D2	Y0590026	≤ -70°C	31 DEC 2027
S-PLEX Enhance E3	Y0570027	≤ -70°C	31 DEC 2027

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator	
Spot Number	Name	Recovery	CV
1	Tau (pT217)	80% - 120%	< 20%

		LLOQ				LLOD	
Name	Range Tested	Observed	Units	Recovery	CV	Observed	Units
Tau (pT217)	1050 - 16800	1050	fg/mL	80% - 120%	< 20%	304	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
Tau (pT217)	20X Calibrator	74600000	fg/mL

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 SEP 2025